

# Sartocon® 2 Plus Stainless Steel Holder

Tangential Flow Filtration  
Filter Holder for  
Sartocon® Cassettes



## Description

### Sartocon® 2 Plus

The Sartocon® 2 Plus stainless steel holder is optimized for the use with up to ten Sartocon® production scale tangential flow filtration cassettes or two Sartocubes® (max. 7 m<sup>2</sup>). It is designed for applications from 30 L to 1,000 L volume.

### Target use:

- Process development
- Preclinical trials
- Clinical trials
- Pilot lots
- Production

Sartocon® cassettes with a filter area of up to 0.7 m<sup>2</sup> each, are available with Polyethersulfone and Hydrosart® membranes.

The Sartocon® 2 Plus system covers the whole range of pharmaceutical and biotechnological Tangential Flow Filtration (TFF) applications like concentration and diafiltration of proteins, vaccines, viruses, antibodies, oligo nucleotides, endotoxin removal etc. The system is also suitable for cell harvesting and clarification processes. The family of Sartorius Stedim Biotech TFF filter holders feature the latest advances in TFF design:

- Vertical orientation of cassettes
- Self draining systems
- Autoclavable
- No welds
- Minimized hold up volume

The Sartocor® 2 Plus holding device is part of a modular micro- and ultrafiltration system and fits up to ten Sartocor® filter cassettes. It is designed for process development and pilot scale production in pharmaceutical and biotechnology applications. Easy handling design and high quality materials make this holder the perfect tool.

The Sartocor® cassettes are placed between filter and clamping plate without a need for additional gaskets. It's not even necessary to remove the clamping plate for installation of the cassettes. Just place the cassettes on the guide rods and close the holder with a torque wrench.

The vertical positioning of the cassettes allows complete draining of retentate and permeate channels by gravity during harvesting procedures.

## Features and Benefits

Feature	Benefits
Cassettes in vertical position	Self draining
Tri-clamp connectors	Safe fit of accessories
No welds	No corrosion
Fits up to ten Sartocor® cassettes	High flexibility for up to 1,000 L working volume

The Sartocor® 2 Plus's flow distribution plate is machined from a single piece of stainless steel. It's sanitary 1 ½" tri-clamp connections are not welded but machined from the same plate. This eliminates potentially problems associated with welds and ensures long service life and safe and reliable connection to other equipment.

The feed and retentate connectors are located on one side of the plate. This allows easy integration into any skid design and assures a compact system with low minimum working volume. Connectors are situated to eliminate potential air locks and to assure thorough cleaning and total product recovery. The Sartocor® 2 Plus device is available a mechanical polished version.

## Technical Specifications

Materials of Construction	
Filter and Clamping plate	316 L stainless steel
Connectors	1" - 1 ½" sanitary Tri-Clamp flange (DN25)
Nuts for tie rods	Nickel coated bronze
Other components	316 L Stainless steel

### Technical Data

Surface finish	Ra ≤ 0.6 µm
Maximum number of cassettes	10 Sartocor® cassettes
Dimensions (L × W × H)	371 × 270 × 348 mm
Weight	Approx. 50 kg

# Ordering Informations and Accessories

Sartocon® 2 Plus microfiltration set	17546---201
Sartocon® 2 Plus ultrafiltration set	17546---202
Torque wrench for Sartocon® 2 Plus holder	17128
Socket, 27 mm	6986135
Pressure gauge, 0-6 bar, 1-1 ½" autoclavable	17546---003
Diaphragm valve, 1-1 ½"	17546---005
Silicone gasket (FDA), 1-1 ½"	17546---012
Clamp for 1-1 ½" Tri-Clamp	SB-18-00-0194
Sartocon® 2 Plus manifold for permeate outlet	17546---016
Hose connector to 1-1 ½" Tri-Clamp flange	17546---018
Sartocon® Adapterplate (round to oval adapter)	302020A-C
Sartocon® Single Use Adapter Plates	17546---113-C
Sartocon® Flush Plate PP-DWST	302030C-C
Sartocon® Dummy Modules PP-DWST	
20 mm	302020D-C
40 mm	302040D-C
100 mm	302100D-C

## Sartocon® 2 Plus, AISI 316 L (1.4435)

Order number	Standard	F   R (mm)	P (mm)	Mirror-inverted
17546DD260160E	DIN	26	16	
17546DZ221157E	Zoll	22.1	15.7	
17546DI297181E	ISO	29.7	18.1	
17546DJ230230E	JIS-G	23	23	
17546DD260160EM	DIN	26	16	■
17546DZ221157EM	Zoll	22.1	15.7	■
17546DI297181EM	ISO	29.7	18.1	■
17546DJ230230EM	JIS-G	23	23	■

F - Feed  
R - Retentate  
P - Permeate

**Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

**USA**

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

 **For further contacts, visit**  
[www.sartorius.com](http://www.sartorius.com)